



Renewable Electricity Certificate

British Gas Trading Ltd certifies that 100% of electricity supplied by British Gas to

Promega UK Limited

will be backed by **Renewable Energy Guarantees of Origin** for the period:

From:

1 October 2024

To:

30 September 2026



Matt Wood
Director, British Gas Business



SALE/GEC/002

* 100% of the electricity supplied is renewable, backed by certificates of renewable energy guarantees of origin. The emission factor conforms with the GHG Protocol Scope 2 guidance and may be used for the 'market-based' method.